

Labeling of *in situ* hybridization DNA probes

- > Fluorescent dUTP & dCTP
- > Hapten-modified dUTP & dCTP
- > CLICKable dUTP & dCTP

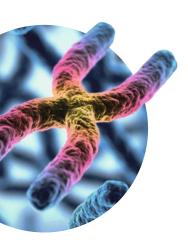
ATAAGC AACO TTCAGGGAAGAA CTAUAACTGCCAC

> Amine-modified dUTP & dCTP



IFTA AG Certified QMS and EMS according to DIN EN ISO 9001 and DIN EN ISO 14001 Reg.-No.: ICV03597 034 and ICV03597 534

Probes & Epigenetics



Introduction

DNA fragments of variable length ("DNA probes") are able to form non-covalent, highly specific duplexes with a complementary nucleic acid strand (termed hybridization). When a label is attached to such a hybridization probe it can efficiently serve for the detection of a defined DNA target sequence.

Labeled DNA probes are routinely used in research and clinical diagnostics for *in situ* hybridization experiments where specific chromosomal DNA sequences as well as their potential aberrations (mutations, deletions and/or duplications) are detected and localized within fixed tissues and cells.

Visualization of DNA *in situ* is either performed by a fluorescence read-out (Fluorescence *in situ* hybridization (FISH)) or by chromogenic detection (Chromogenic *in situ* hybridization (CISH)) that relies on an enzymatic reaction causing precipitation of colored substrates.

Labeling of DNA probes for *in situ* hybridization is most often performed by enzymatic incorporation of labeled dNTPs using *Taq* DNA polymerase (= Polymerase Chain Reaction (PCR)) or a mixture of DNase I/DNA polymerase I (=Nick Translation) (Tab. 1) ^[1,2].

Table 1 DNA probes for in situ hybridization can be enzymatically labeled via Nick Translation or PCR.

	Nick Translation	PCR
Principle	eIntroduction of random single-strand breaks ("nicks") into dsDNA by DNAse I that are subsequently filled by DNA Polymerase I using a labeled dNTP as substitute for its natural counterpart.Amplifi (PCR w loci (PC primer incorp.	
Template	pmol amounts of linearized dsDNA	fmol amounts of linearized dsDNA
Size of labeled fragment	100–500 bp dsDNA	up to 1500 bp
DNA amplification	no	yes

The most prominent labeled dNTP is dUTP labeled at the C5 position (Fig. 1) which is incorporated as partial substitute for its natural counterpart dTTP. Depending on the template sequence (e.g GC-rich) however, the usage of labeled dCTP or a combination of both might be favorable for achieving the desired labeling.

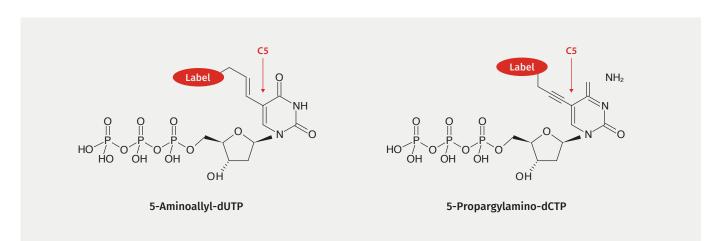


Figure 1 dUTP and dCTP labeled at the C5 position via aliphatic aminoallyl or propargylamino linker, respectively, are compatible with enzymatic incorporation, show minimized influence on Watson-Crick base pairing and thus yield good hybridization results. Label: fluorescent dye, hapten (e.g. Biotin, Digoxigenin, DNP) or reactive group such as amine or CLICKable moiety.

Probe labeling can be performed in a one-step or two step procedure (Fig. 2). The labeling strategy depends on the required sensitivity for the downstream application and the suitability of an appropriately labeled dUTP or dCTP for an enzymatic incorporation method.

For example, labeling with near-infrared dyes is of great importance when working with biological samples since near IRfluorescence is not disturbed by biological autofluorescence (haemoglobins or cytochromes). Nucleotides labeled with near IR-dyes however, often show insufficient labeling efficiencies caused by the bulky fluorophore. Furthermore, dUTP or dCTP labeled with hydrophobic fluorescent dyes often show poor PCR incorporation efficiencies compared to hydrophilic dyes.

These obstacles can be circumvented by a) use of smaller labels such as haptens (Fig. 2A) or b) use of dUTPs/dCTPs carrying a reactive group that is coupled with the desired fluorescent dye in a second step (Fig. 2B).

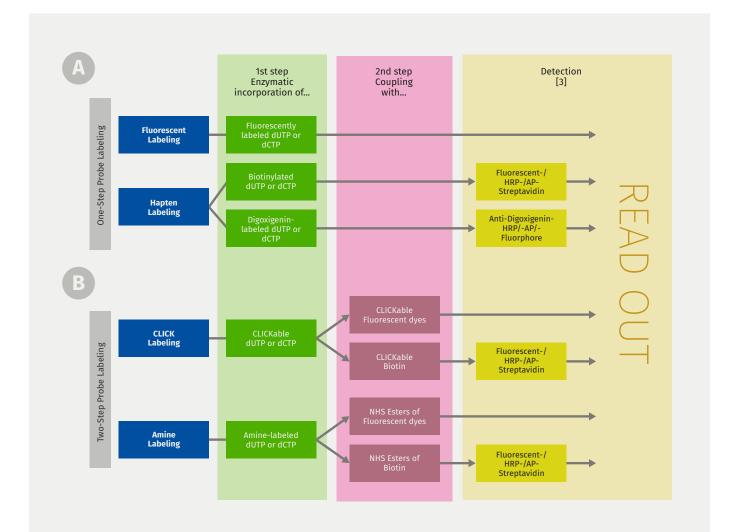


Figure 2 One-step probe labeling (A) is performed by incorporation of dUTP or dCTP carrying a label (dye or hapten) while two-step probe labeling (B) requires first introduction of a reactive group (amine or "Clickable" moiety) followed by coupling with an appropriate label. HRP: horse radish peroxidase, AP: alkaline phosphatase, NHS: N-hydroxy-succinimide.

Recommended Reading:

[1] Morrison et al. (2003) Labeling Fluorescence In situ Hybridization Probes for Genomic Targets. Methods in Molecular Biology **204**: 21.

[2] Wiegant et al. (1997) Probe Labeling and Fluorescence In situ Hybridization. In: Current Protocols in Cytometry 8.3.1, John Wiley & Sons Inc.

BUY ONLINE →

Fluorescent dUTP & dCTP

Fluorescent dUTP and dCTP can be used for enzymatic labeling of DNA probes in a one-step procedure (Fig. 2). The labeled DNA probe is directly visualized by fluorescence microscopy^[3] or spectroscopy. Combinatorial labeling, e.g. labeling of probes with different or a unique set of fluorophores, allows simultaneous visualization of multiple targets (multiplexing)^[4–7].

The choice of fluorophore(s) depends on the available excitation source, filter sets, application (single-colour or multi-colour imaging) (Tab. 2) and desired enzymatic labeling technique (Nick Translation or PCR) (Tab. 3). Both photostability and hydrophilicity of a fluorophore are critical parameters that should be taken into account during decision making. For example, fluorophores with improved photostability may generally increase sensitivity thus improving detection limits of target sequences. The hydrophilicity of a fluorophore however, directly affects the substrate properties of correspondingly labeled dUTP & dCTP.

dUTP & dCTP labeled with hydrophilic (e.g. Cy3) or moderately hydrophobic dyes (e.g. Texas Red) are ideally suited for both PCR and Nick Translation labeling. In contrast, hydrophobic fluorophores (e.g. ATTO647N) cause an early termination of DNA amplification, most likely due to dye-dye or dye-enzyme interactions, and are thus recommended for Nick Translation only.

Table 2 Spectral properties of selected fluorescent dyes. ++: good, +: moderate, -: poor.

Please note: The choice of fluorophore depends on the final application and is a compromise between hydrophilicity and photostability that determine enzymatic incorporation efficiency and detection sensitivity, respectively.

Emission colour	Буе	Exc _{max} [nm]*	Em _{max} [nm]*	ε _{max} [L × mmol⁻¹ cm⁻1]**	Substitute for	Hydrophilicity	Photostability
	AMCA	350	450	19.0		+	+
Blue	DEAC	426	480	57.0		-	+
	ATTO425	436	484	45.0	DEAC	+	++
	Rhodamine	505	530	85.0		_	+
	Fluorescein	492	517	83.0		++	-
Green	ATTO488	501	523	90.0	Fluorescein, Rhodamine, Alexa Fluor 488, Cy2	++	++ (>fluorescein, rhodamine, Cy2)
	AF488	494	515	73.0	Fluorescein, Rhodamine, ATTO488, Cy2	++	++ (>fluorescein, rhodamine, Cy2)
Yellow-Green	ATTO532	532	553	115.0	Alexa Fluor 532	++	++
	Cy3	550	570	150.0		++	+
Yellow	ATTO550	554	576	120.0	Cy3	_	++ (> Cy3)
Overage	Texas Red	588	609	80.0	Су3.5	+	+
Orange	AF594	590	617	92.0	Texas Red	++	++
	J647	637	670	94.0	Cy5, ATTO647N, Alexa Fluor 647	++	++ (> Cy5)
Red	ATTO647N	644	669	150.0	Cy5, Alexa Fluor 647	_	++ (> Cy5)
	Cy5	649	670	250.0		++	_

4

5

Table 3 Enzymatically incorporable fluorescent dUTP and dCTP. n/a: not applicable

		Recommended dTTP or dCTP substitution by modified dNTP [%]*					
Nucleotide	Cat. No.	PCR (Taq Pol)	Nick Translation (DNAse I , DNA Pol I)				
dUTP							
AMCA-6-dUTP NU-803-AMCA 20-30% 30-50%							
DEAC-dUTP	NU-803-DEAC	n/a	30-50 %				
dUTP-ATTO425	NU-803-425	20-30 %	30-50 %				
Rhodamine-12-dUTP	NU-803-RHOX	n/a	30-50 %				
Fluorescein-12-dUTP	NU-803-FAMX	30-50 %	30-50 %				
dUTP-XX-ATTO488	NU-803-XX-488	30-50 %	30-50 %				
dUTP-XX-ATTO532	NU-803-XX-532	30-50 %	30-50 %				
dUTP-Cy3	NU-803-Cy3	30-50 %	30-50 %				
dUTP-ATTO550	NU-803-550	n/a	30-50 %				
dUTP-Texas Red	NU-803-TXR	20-30 %	30-50 %				
dUTP-J647	NU-803-J647	30-50 %	30-50 %				
dUTP-Cy5	NU-803-Cy5	30-50 %	30-50 %				
dUTP-ATTO647N	NU-803-647N	n/a	30-50 %				
	d	СТР					
Fluorescein-12-dCTP	NU-809-FAMX	30-50 %	30-50 %				
Rhodamine-12-dCTP	NU-809-RHOX	n/a	30-50 %				
dCTP-Cy3	NU-809-Cy3	30-50 %	30-50 %				
dCTP-ATTO550	NU-809-550	n/a	30-50 %				
dCTP-Texas Red	NU-809-TXR	20-30 %	30-50 %				
dCTP-J647	NU-809-J647	30-50 %	30-50 %				
dCTP-Cy5	NU-809-Cy5	30-50 %	30-50 %				
dCTP-ATTO647N	NU-809-647N	n/a	30-50 %				

*Please note:

Optimal final concentration of the dye-labeled dNTP may vary depending on the application and assay conditions. For optimal product yields and optimal incorporation rates an individual optimization of the dye-labeled dUTP/ dTTP or dye-labeled dCTP/dCTP ratio is recommended.

dTTP is substituted by modified dUTP, dCTP is substituted by modified dCTP

Recommended reading:

[3] Jena Bioscience GmbH (a): Non-radioactive Labeling of DNA and RNA, http://www.jenabioscience.com/images/741d0cd7d0/bro_FluorescentProbes_WEB.pdf, 23.04.2015.

- [4] Speicher et al. (1996) Karyotyping human chromosomes by combinatorial Multi-Fluor FISH. Nat. Genetics 12:368.
- [5] Fauth et al. 2001) Classifying by colors: FISH-based genome analysis. Cytogen. Cell Genet. 93:1.

[6] Bayani et al. (2004) Multi-color FISH techniques. Curr. Protoc. Cell. Biol. 22: Unit 22.5.

[7] Anderson et al. (2004) Multiplex fluorescence in situ hybridization (M-FISH). Methods Mol. Biol. 659:83.

Website fluorescent PCR Labeling Kits at ↗



Website fluorescent Nick Translation Labeling Kits at 🔻



6

BUY ONLINE →

Hapten-modified dUTP & dCTP

Hapten-modified dUTP & dCTP are suitable for enzymatic labeling of DNA probes in a one-step procedure (Fig. 2) using reporter molecules for subsequent detection^[8].

Such reporter molecules are either coupled with a fluorophore for visualization by fluorescence microscopy (Fluorescence *in situ* hybridization (FISH)) or with a reporter enzyme that generates signals through an enzymatic reaction with chromogenic substrates (Chromogenic *in situ* hybridization (CISH)) with subsequent detection by bright-field microscopy^[8]. This signal amplification usually results in higher sensitivity of indirect detection methods compared to direct fluorescent labeling.

The most commonly used haptens are Biotin, Digoxigenin and Dinitrophenol (DNP). Accordingly labeled dUTP & dCTP are excel-

lent enzymatic substrates (Tab. 4)^[9-12] whose incorporation and detection efficiency by secondary reporter molecules is strongly influenced by linker length.

In general short linkers result in high probe yields but may prevent sufficient accessibility of label on the probe surface while long linkers result in increased label accessibility towards detection reagents but result in lower probe yields. An optimal balance between incorporation rate and detection efficiency is achieved by an 11- atom or 16-atom linker^[12].

For most applications Biotin is the preferred label however, in some cases endogenous Biotin from biological samples may interfere with detection, resulting in high backgrounds or even false positives. In such cases, the use of Digoxigenin, a steroid exclusively present in Digitalis plants, is recommended.

Table 4 Enzymatically incorporable Hapten-modified dUTP and dCTP. n/a: not applicable

		Recommended dTTP or dCTP substitution by modified dNTP [%]*		
Nucleotide	Cat. No.	PCR (Taq Pol)	Nick Translation (DNAse I , DNA Pol I)	
	dUJ	P		
Biotin-11-dUTP	NU-803-BIOX	30-50 %	30-50 %	
Biotin-16-dUTP	NU-803-BIO16	30-50 %	30-50 %	
Digoxigenin-11-dUTP	NU-803-DIGX	30-50 %	30-50 %	
DNP-dUTP	NU-803-DNP	n/a	30-50 %	
	dCT	Р		
Biotin-11-dCTP	NU-809-BIOX	30-50 %	30-50 %	
Biotin-16-dCTP	NU-809-BIO16	30-50 %	30-50 %	

*Please note:

Optimal final concentration of the Hapten-labeled dNTP may vary depending on the application and assay conditions. For optimal product yields and optimal incorporation rates an individual optimization of the Hapten-labeled dUTP/ dTTP or Hapten-labeled dCTP/dCTP ratio is recommended.

dTTP is substituted by modified dUTP, dCTP is substituted by modified dCTP

Recommended reading:

- [8] Jena Bioscience GmbH (a): Non-radioactive Labeling of DNA and RNA, http://www.jenabioscience.com/images/741d0cd7d0/bro_FluorescentProbes_WEB.pdf, 23.04.2015
- [9] Anderson et al. (2005) Incorporation of reporter-labeled nucleotides by DNA polymerases. Biotechniques 38:257.

[10] Jackson *et al.* (1991) Detection of Shiga Toxin-Producing Shigella dysenteriae Type 1 and Escherichia coli by Using Polymerase Chain Reaction with Incorporation of Digoxigenin-11-dUTP. J Clin Microbiol. **29(9)**:1910.

[11] Dauwerse et al. (1999) Two-colour FISH detection of the inv(16) in interphase nuclei of patients with acute myeloid leukemia. Br J Haematol 106:111.

[12] Jena Bioscience GmbH (b): Biotin and Digoxigenin Labeled Nucleotides, http://www.jenabioscience.com/images/741d0cd7d0/Biotin_DIG_Nuleotides.pdf, 23.04.2015.

Website Biotin and Digoxigenin PCR Labeling Kits at ↗



Website Biotin and Digoxigenin Nick Translation Labeling Kits at ↗



CLICKable dUTP & dCTP

CLICKable dUTP & dCTP (Tab. 5) can be used for both hapten and fluorescent DNA labeling by a two-step procedure (Fig. 3).

The small-sized CLICK reactive moieties (DBCO or Azide) possess excellent substrate properties resulting in incorporation efficiencies of CLICKable dUTP & dCTP analogous to their Amine-modified counterpart (Tab. 5). CLICK labeling however, offers a superior labeling procedure compared to strongly pH-dependent Amine-NHS-Ester-based reactions that are often hampered by competing NHS ester hydrolysis.

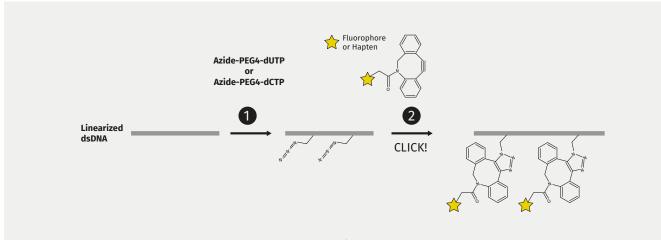


Figure 3 CLICK DNA Labeling is performed by a two-step procedure. 1) Enzymatic Azide-functionalization of dsDNA by PCR or Nick Translation. 2) Copper-free CLICK-labeling via a DBCO-functionalized detection molecule. DBCO-PEG4-dUTP and –dCTP are similarly incorporated but labeled with Azide-functionalized detection molecules.

Table 5 Enzymatically incorporable CLICKable dUTP & dCTP.

		Recommended dTTP or dCTP substitution by modified dNTP [%]*				
Nucleotide	Cat. No.	PCR (Taq Pol)	Nick Translation (DNAse I , DNA Pol I)			
	dUTP					
Azide-PEG4-dUTP	NU-1705	30-50 %	30-50 %			
DBCO-PEG4-dUTP	CLK-T09	30-50 %	30-50 %			
	dCTP					
Azide-PEG4-dCTP	CLK- 061	30-50 %	30-50 %			
DBCO-PEG4-dCTP	CLK- 060	30-50 %	30-50 %			

*Please note:

Optimal final concentration of the CLICKable dNTP may vary depending on the application and assay conditions. For optimal product yields and optimal incorporation rates an individual optimization of the CLICKable dUTP/ dTTP ratio is recommended.

dTTP is substituted by modified dUTP, dCTP is substituted by modified dCTP

Recommended reading:

[13] Jena Bioscience GmbH: CLICK Chemistry Background Information, http://www.jenabioscience.com/images/741d0cd7d0/20140306_Click_Chemistry_Background_information_pdf_creator. pdf, 06.01.2016.

Website Azide and DBCO detection molecules at ↗



Website CLICK PCR Labeling Kits at ↗



Website CLICK Nick Translation Labeling Kits at ↗



BUY ONLINE →

Amine-modified dUTP & dCTP

Amine-modified dUTP & dCTP are historically used for both hapten and fluorescent DNA labeling by a two-step procedure (Fig. 4).

Step 1 consists of the enzymatic incorporation of the Aminemodified dNTP (Tab. 6) which is subsequently labeled with an Amine-reactive (e.g. NHS ester) fluorescent dye or hapten^[14–15].

Table 6 Enzymatically incorporable Amine-modified dUTP and dCTP.

Due to the small size of the amino group, higher incorporation efficiencies may be achieved compared to enzymatic labeling with fluorescent dUTPs and dCTPs – especially when carrying bulky or hydrophobic fluorescent dyes.

		Recommended dTTP or dCTP substitution by modified dNTP [%]*			
Nucleotide	Cat. No.	PCR (Taq Pol) Nick Translation (DNAse I , DNA Pol			
dutp					
Aminoallyl-dUTP	NU-803	30-50 %	30-50 %		
dCTP					
Propargylamino-dCTP	NU-809	30-50 %	30-50 %		

*Please note:

Optimal final concentration of the Amine-labeled dNTP may vary depending on the application and assay conditions. For optimal product yields and optimal incorporation rates an individual optimization of the Amine-labeled dUTP/ dTTP or Amine-labeled dCTP/dCTP ratio is recommended. dTTP is substituted by modified dUTP, dCTP is substituted by modified dCTP

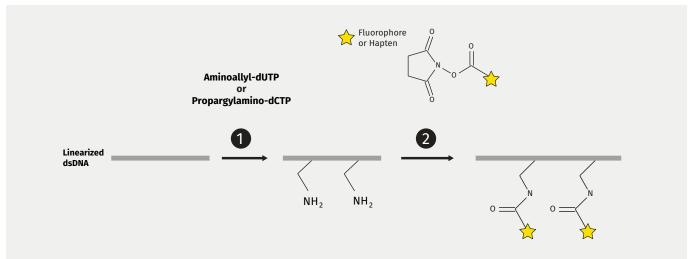


Figure 4 Amine DNA Labeling is performed by a two-step procedure. 1) Enzymatic Amine-functionalization of dsDNA by PCR or Nick Translation. 2) Amine Labeling via a N-Hydroxysuccinimid (NHS)-functionalized detection molecule.

Recommended reading:

[14] Dirsch *et al.* (2007) Probe production for *in situ* hybridization by PCR and subsequent covalent labeling with fluorescent dyes. *Appl. Immunohistochem. Mol. Morphol.* 3:332.
 [15] Cox *et al.* (2004) Fluorescent DNA hybridization probe preparation using amine modification and reactive dye coupling. *BioTechniques* 36:114.

Website NHS Ester detection molecules at ↗



8

BUY ONLINE -

Jena Bioscience is a primary nucleotide manufacturer offering bulk amounts as well as custom formulations, packaging & labeling.

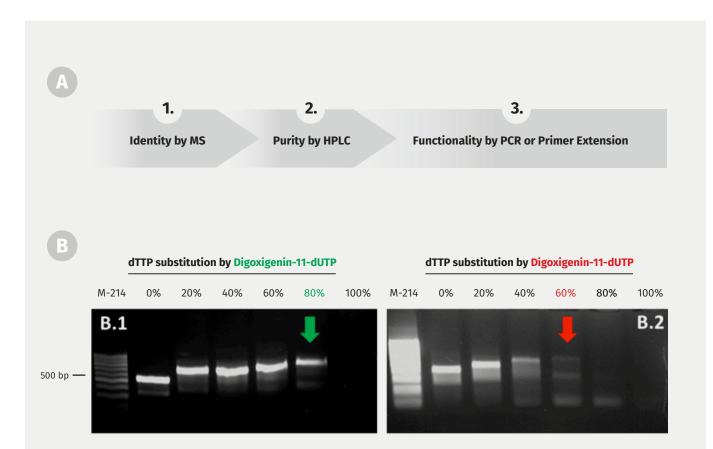


Figure 5 Quality control of selected labeled nucleotides. A) Three-step quality control procedure B) Functional Quality Control Example (Step 3): Each lot of Digoxigenin-11-dUTP is – besides a number of other tests – analyzed in a PCR reaction with increasing substitution of dTTP by Digoxigenin-11-dUTP (0–100%). The band shift indicates Digoxigenin-11-dUTP incorporation. Lots yielding a visible PCR prod-uct at 80% substitution of dTTP by Digoxigenin-11-dUTP by Digoxigenin-11-dUTP are accepted (B.1). B.2: Non-Conforming Digoxigenin-11-dUTP lot (source: other supplier).



IFTA AG Certified QMS and EMS according to DIN EN ISO 9001 and DIN EN ISO 14001 Reg.-No.: ICV03597 034 and ICV03597 534 Jena Bioscience has established a quality management system according to DIN EN ISO 9001 and an environmental management system according to DIN EN ISO 14001. We commit ourselves to international standards and continuously improve our quality and environmental management system and our inhouse processes to meet our high quality standards.

Terms and Conditions of Sales

Ordering

You have the following options when ordering products directly from Jena Bioscience:

- Mail orders
- \cdot Telephone orders
- \cdot 24 hour fax ordering
- Online ordering

Please provide us with the following information when ordering:

- \cdot Your name, name of institution
- $\boldsymbol{\cdot}$ Billing and shipping address
- PO number (if applicable)
- Catalog number of products and quantities needed
- Contact person and contact data for questions

Mail orders

Please send your mail orders to the following address:

Jena Bioscience GmbH Loebstedter Strasse 71 07749 Jena, Germany

Telephone orders

We will accept telephone orders from Monday to Friday between 8:00 am and 16:00 pm Central European Time. +49 - 3641 - 628 5000

24 hour fax ordering

Please send your fax order to: +49 - 3641 - 628 5100

Online ordering

Jena Bioscience products can be ordered online. When ordering by e-mail, please direct your orders to orders@jenabioscience.com

Products can also be ordered online in our online shop: http://www.jenabioscience.com

Important Notice:

Products that have been ordered by mistake cannot be returned to Jena Bioscience. Products that are returned unrequestedly to Jena Bioscience will not be accepted, but fully charged to the customer's account.

Shipping

The standard shipping term is CPT (Incoterms 2010). Jena Bioscience decides on the shipping carrier.

Alternatively, customer may arrange shipping himself or elect a carrier of his choice, e.g. by providing his carrier account number.

In this case, products will be shipped at customer's risk and Jena Bioscience will not be liable for loss, damage or thawing of products.

Prices and Charges

Please note that the prices of products in the catalog and on our website do not include freight charges, duties, taxes or customs fees.

Freight charges will be prepaid and added to the invoice. Freight charges for online orders are indicated when you check out of the online store. If you need information on freight charges for your particular order, please contact us with all necessary information.

Jena Bioscience will not pay any duties, taxes or customs fees.

Products and prices are subject to change without notice. Current pricing will be confirmed at the time of your order. No minimum order required.

We offer free shipping of all orders worth EUR 250+ (excl. VAT) within Germany.

Payment

Invoices will be issued after your order has been shipped and will be sent to the billing address by separate mail. Invoices will not be included within the shipments. In case of partial deliveries, separate invoices will be issued after each shipment has left Jena Bioscience. You will find payment information (bank addresses and account data) on each invoice. Jena Bioscience accepts payment by:

Check

Please send your payment checks to the following address: Jena Bioscience GmbH Loebstedter Strasse 71 07749 Jena, Germany

We kindly ask you to make sure that our invoice number and your customer number appear on the check.

Wire transfer

All information necessary for wire transfers will be shown on each invoice.

Our VAT number (for EU customers): DE 195825742

Credit card

Jena Bioscience accepts the following credit cards:

- VISA
- Mastercard
- American Express

If you wish to pay by credit card, please provide the following credit card information:

- \cdot Card holder
- Card number
- \cdot Expiry date
- Security code (VISA / Mastercard: 3 digits, to be found on your card's back side in the upper right corner of the signature field; AmEx: usually 4 digits (sometimes only three), to be found on the front side of your card above the card number)

Patent Disclaimer

Unless explicitly stated, no license or immunity under any patent is either granted or implied by the sale of any of our products. Jena Bioscience does not warrant that the resale or use of its products delivered will not infringe the claims of any patent, trademark or copyright covering the use of the product itself or its use in the operation of any process. Furthermore, the purchaser assumes all risks of patent, trademark or copyright infringement associated with any such use, combination or operation.

11

Billing address	Shipping address
Customer number	Name
University/Company	University/Company
Institute/Department	Institute/Department
Address	Address
Postcode	Postcode
City/Signature	City/Signature
VAT number (EEC only)	Phone
PO number	Fax
Date / Signature	Email

If you wish to pay by credit card, please provide the following credit card information: I want to pay by



```
Card holder
```

Card number

Expiry date

Security code

(VISA / Mastercard: 3 digits on card's back side, upper right corner of signature field; AmEx: 4 digits, card's front side, above card number)

	Catalog number	Product	Quantity	Net Price per Item EURO	Net Price all Items EURO
1					
2					
3					
4					
5					
6					
7					
8					
9					
10					
11					
12					
13					
14					
	1		1	Total	



Jena Bioscience GmbH Loebstedter Str. 71 07749 Jena Germany Phone +49(0)3641-6285 000 Fax +49(0)3641-6285 100 info@jenabioscience.com www.jenabioscience.com



Jena Bioscience GmbH Loebstedter Str. 71 07749 Jena Germany

Phone +49(0)3641-6285000 Fax +49(0)3641-6285100 info@jenabioscience.com

www.jenabioscience.com



IFTA AG Certified QMS and EMS according to DIN EN ISO 9001 and DIN EN ISO 14001 Reg.-No.: ICV03597 034 and ICV03597 534

CCACGAAAGGGAA ATAAGC AACO TTCAGGGAAGAA CTAUAACTGCCAC ACCCACGAAAGGGAA ATAAGC AACO TTCAGGGAAGAA CAGGGAAGAA CTAUAACTGCCAC **ACCCACGAAAGGGAA ATAAGC AACO TTCAGGGAAGAA CTAUAACTGCCAC** ACCCACGAAAGG AAGGGAA ATAAGC AACO TTCAGGGAAGAA CTAUAACTGCCAC ACCCACGAAAGGGAA ATAAGC AACO TTCAGGGAAGAA CTAUA

Probes & Epigenetics